

PE Anti-human CD70 Antibody *Ki-24*Catalog number: 107001M0, 107001M1, 107001M2
Unit size: 25 tests, 100 tests, 500 tests**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG3 kappa
Immunogen	CD70 (CD27 ligand, CD27L, TNFSF7)
Clone	Ki-24
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE
Excitation Wavelength	566 nm
Emission Wavelength	574 nm

Applications

The Ki-24 monoclonal antibody binds to human CD70, a 50 kD member of the tumor necrosis factor superfamily frequently found on the surface of T cells, dendritic cells, B cells and natural killer cells. CD70 is associated with a variety of biologically interesting macromolecules/ligands, for instance, CD27 ligand. CD70 is a fairly uncommon antibody target, with a little more than 2900 publications in the last decade. Even still, CD70 has been widely used in immunology and costimulatory molecules research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE (ex/em = 566/574 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).